

Catalog #	Aliquot Size
<b>M821-40N-20</b>	<b>20 µg</b>
<b>M821-40N-100</b>	<b>100 µg</b>
<b>M821-40N-1000</b>	<b>1 mg</b>

## Recombinant Human CCL2

Recombinant human protein expressed in E. coli cells

**Catalog # M821-40N**

Lot # G3211-11

### Product Description

Recombinant Human CCL2 was expressed in E. coli cells. The protein accession number is [P13500](#)

### Gene Aliases

Monocyte Chemoattractant Protein-1, MCP-1, CCL2, MCAF, C-C motif chemokine 2, HC11, Monocyte chemoattractant and activating factor, Monocyte secretory protein JE, Small-inducible cytokine A2, MCP1, SCYA2, Recombinant Human Monocyte Chemoattractant Protein-1 (CCL2)

### Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

### Formulation

MCP-1 was lyophilized from PBS pH 7.4.

### Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

### Storage and Stability

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

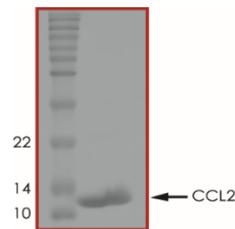
### Scientific Background

MCP-1 belongs to the family of chemotactic cytokines. MCP-1 is expressed by monocytes, vascular endothelial cells, smooth muscle cells, glomerular mesangial cell, and osteoblastic cells. MCP-1 has been shown to exhibit biological activities other than chemotaxis. It can induce the proliferation and activation of killer cells known as CHAK. Elevated levels of MCP-1 are observed in atherosclerotic plaques that are rich in macrophages. The factor activates the tumoricidal activity of monocytes and macrophages in vivo. It regulates the expression of cell surface antigens (CD11c, CD11b) and the expression of cytokines (IL-1, IL-6). MCP-1 is a potent activator of human basophils, inducing degranulation and the release of histamines. In basophils activated by IL-3, IL-5 or GM-CSF MCP-1 enhances the synthesis of leukotriene C4.

### Specific Activity

Determined by its ability to chemoattract human monocytes using a concentration range of 2-40 ng/mL.

### Purity



Recombinant Human CCL2 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **9.0 kDa**

## Recombinant Human CCL2

Recombinant human protein expressed in E. coli cells

Catalog Number	M821-40N
Specific Lot #	G3211-11
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**